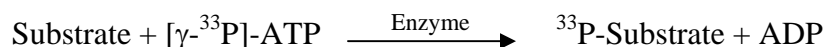


JNK3

(Mitogen-activated protein kinase 10, Stress-activated protein kinase JNK3, c-Jun N-terminal kinase 3, MAP kinase p49 3F12, MAPK10, JNK3A, PRKM10, MK10)

- **CAT#:** JNK3
- **Enzyme:** Human JNK3
- **Substrate:** ATF2, 3 μ M
ATP 10 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	35,000
JNK Inhibitor VIII	6,900

- **Scientific Information:** P53779 (<http://www.expasy.org/uniprot/P53779>)

Reference compound IC50 for JNK3

